

# CMP-Neu5Ac

Nucleotide sugar and the donor substrate for sialyltransferases (SiaTs)

BLG-C408-50 (5 x 10  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

## Ordering Information

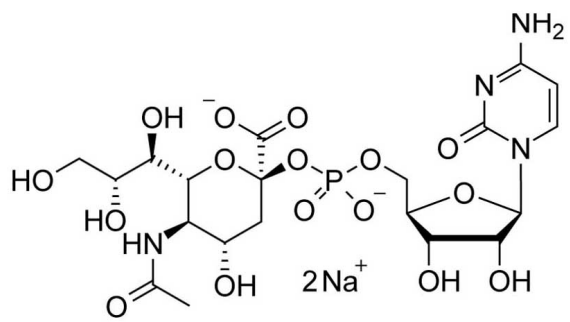
[Order Online »](#)

BLG-C408-10	10 $\mu$ mol
BLG-C408-50	5x10 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)

CMP- $\beta$ -D-N-Acetylneuraminic acid (CMP-Neu5Ac) is a nucleotide sugar and the donor substrate for eukaryotic and prokaryotic sialyltransferases. Furthermore CMP-Neu5Ac is the essential substrate for sialyltransferases (SiaTs) in the synthesis of glycoconjugates.



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** Cytidine-5'-O-monophospho-N-acetylneuraminic acid . sodium salt, CMP-NAN, CMP-NANA, CMP-Sialic acid

**CAS** 3063-71-6

**Formula**  $C_{20}H_{31}N_4O_{16}P$  (free acid)

**MW** 614.5 (free acid)

**Purity** > 95% HPLC

**Quantity** 10 µmol are equivalent to ~ 6.2 mg



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)